

# PRODUCT INFORMATION

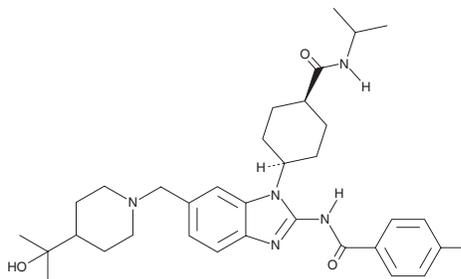


## Belizatinib

Item No. 28458

**CAS Registry No.:** 1357920-84-3  
**Formal Name:** 4-fluoro-N-[6-[[4-(1-hydroxy-1-methylethyl)-1-piperidiny]methyl]-1-[cis-4-[[[(1-methylethyl)amino]carbony]cyclohexyl]-1H-benzimidazol-2-yl]-benzamide

**Synonym:** TSR-011  
**MF:** C<sub>33</sub>H<sub>44</sub>FN<sub>5</sub>O<sub>3</sub>  
**FW:** 577.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 324 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Belizatinib is supplied as a crystalline solid. A stock solution may be made by dissolving the belizatinib in the solvent of choice, which should be purged with an inert gas. Belizatinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of belizatinib in these solvents is approximately 30 mg/ml.

### Description

Belizatinib is an inhibitor of anaplastic lymphoma kinase (ALK) and tropomyosin-related kinases A/B/C (IC<sub>50</sub>s = 0.7 and <3 nM, respectively).<sup>1</sup>

### Reference

1. Sachdev, J., Arkenau, H.T., Infante, J.R., *et al.* Phase (Ph) 1/2a study of TSR-011, a potent inhibitor of ALK and TRK, in advanced solid tumors including crizotinib-resistant ALK positive non-small cell lung cancer. *Eur. J. Cancer* **50(6)**, 165 (2014).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/05/2022

### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM